



PvuI (7)
SgfI (6)
MfeI (82)

1 GGATCTGCATCGCTCCGGTGCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA

101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTGTCTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)

Psp1406I (203)
PvuII (239)
Bsu36I (291)

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCAGCTTTGCTGACCCTGCTTGTCAACTCTACGCTTTTGTTCGTTT

AgeI (552) BspHI (560)

501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCTACCTGAGATCACCGGTATCATGAATCAAAGTCCATTCTGATTTGCTGCCTTATCTT

XbaI (634) 1 ▶ M N Q T A I L I C C L I F

601 TCTGACTCTAAGTGGCATTCAAGGAGTACCTCTCTCTAGAACTGTACGCTGTACCTGCATCAGCATTAGTAATCAACCTGTTAATCCAAGGCTTTAGAA

13 ▶ L T L S G I Q G V P L S R T V R C T C I S I S N Q P V N P R S L E

701 AAACCTGAAATATTCTGCAAGCAATTTTGTCCACGTGTTGAGATCATTGCTACAATGAAAAAGAGGGTGAAGAGATGCTGAAATCCAGAATCGA

47 ▶ K L E I I P A S Q F C P R V E I I A T M K K K G E K R C L N P E S

XmnI (706) **BbrPI (734)** **MscI (865)**

801 AGGCCATCAAGAATTTACTGAAAGCAGTTAGCAAGGAAAGGTCTAAAAGATCTCCTTAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGG

80 ▶ K A I K N L L K A V S K E R S K R S P •

BglIII (847) NheI (859)

901 ACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGTGAATAAACAAGTT

HpaI (997)

MfeI (1008) **EcoRI (1093)**

1001 AACAAACAACATTCATTATTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTCT

1101 AAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAAT

SapI (1275)

1201 GTGCATTAGCTGTTTGCAGCCTCACCTCTTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGTTTGAAGTACTCTTCTTTATGTTTTAAA

SspI (1332) SwaI (1346)

1301 TGCACTGACCTCCACATTCCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAG

EcoO109I (1407)

1401 ATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCAGGCTTCTAG

1501 CTTTAGTTCCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGT

141 ▶ • N R T Y K L P I L E E I T T K V L K G N M E I L V F C D P A Y D

1601 CAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGATAGGGGTGCCTGACAGCCACAATGGTGTCAA

108 ▶ S I L E R C M G C P S V V R I S R D V E D S Y P H R V A V I T D F

StuI (1771)

1701 GTCCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATC

75 ▶ D K Q G N S V A S G I A I A E A C V T V R G I Y A E I H V A S I I

1801 TCCCCAGTCTTGGTCTGATGGCCGCCCGACATGGTCTTGTGCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACAGCTCCAGATCCT

41 ▶ E G T K T R I A A G V H H K N D E Y L M T I K E T A V E V L E L D Q

BbsI (1917)

XmnI (1913) **AseI (1979)**

1901 GCTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCTATAGTGTGATTATACTATGCGGATATACTATGCGGATGATTAATTGTCAAACACGCGTG

8 ▶ Q S I N F T K M

SacI (2036)

2001 GATGGCGTCTCCAGCTTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGA

SpeI (2134)

2101 GTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCGTGAGTCAAAC

SnaBI (2262)

2201 CGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGG

